NPPB

CI⁻ channel blocker

A potent chloride channel blocker (IC $_{50}$ =0.1-100 μ M depending on channel subtype and assay method). Inhibits cyclooxygenase (COX; IC $_{50}$ =8 μ M). Evidence of mitochondrial uncoupling in mouse peritoneal macrophages, due to protonophore activity.

Citations: 7

View Online »

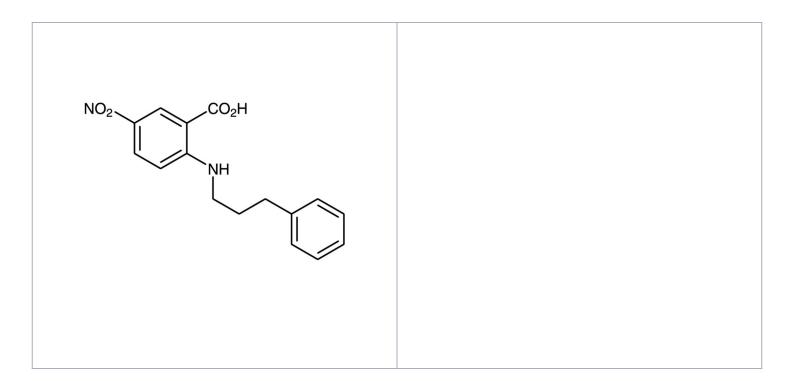
Ordering Information

Order Online »

BML-CM100-0010	10mg
BML-CM100-0050	50mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at room temperature. Stock solutions are

stable for up to 3 months at -20°C.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-Nitro-2-(3'-phenylpropyl-amino)benzoic acid

Appearance Yellow solid.

CAS 107254-86-4

Couple Target Chloride channel, COX

Couple Type Blocker, Inhibitor

Formula $C_{16}H_{16}N_2O_4$

MW 300.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (10mg/ml;

warm).

Last modified: May 29, 2024

